

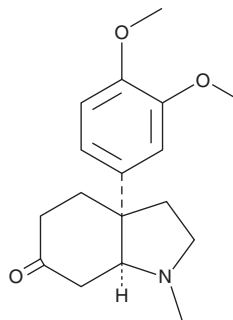
PRODUCT INFORMATION



Mesembrine

Item No. 21100

CAS Registry No.: 24880-43-1
Formal Name: (3aS,7aS)-3a-(3,4-dimethoxyphenyl) octahydro-1-methyl-6H-indol-6-one
Synonyms: Mesembranone, (-)-Mesembrine
MF: C₁₇H₂₃NO₃
FW: 289.4
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Sceletium tortuosum*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Mesembrine is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform purged with an inert gas can be used. The solubility of mesembrine in chloroform is approximately 10 mg/ml.

Description

Mesembrine is an alkaloid that has been found in *S. tortuosum* and has serotonin (5-HT) reuptake inhibitory activity.¹ It selectively binds to the 5-HT transporter (SERT; IC₅₀ = 4.3 µg/ml) over a variety of receptors (IC₅₀s = 148->750 µg/ml).² Mesembrine inhibits the reuptake of 5-HT *in vitro* (K_i = 1.4 nM). It has analgesic effects in the hot plate test when administered at a dose of 20 mg/kg but does not induce a preference in the conditioned place preference test (CPP).^{2,3}

References

1. Krstenansky, J.L. Mesembrine alkaloids: Review of their occurrence, chemistry, and pharmacology. *J. Ethnopharmacol.* **195**, 10-19 (2017).
2. Harvey, A.L., Young, L.C., Viljoen, A.M., *et al.* Pharmacological actions of the South African medicinal and functional food plant *Sceletium tortuosum* and its principal alkaloids. *J. Ethnopharmacol.* **137(3)**, 1124-1129 (2011).
3. Loria, M.J., Ali, Z., Abe, N., *et al.* Effects of *Sceletium tortuosum* in rats. *J. Ethnopharmacol.* **155(1)**, 731-735 (2014).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/23/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM